



CERTIFICATE OF ANALYSIS

Product Description: Human IL-4 Calibrator
Catalog Number: C01A3-2
Lot Number: 590652
Expiration Date: 30 JUN 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-4	133000	fg/mL

Reconstitution Volume: 1000 µL

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		20 AUG 2026